

AlphaLISA®

HEK 293 Host Cell Protein Detection Kit**Product number:** AL3198C **Lot Number:** 3304903**Material provided:****Format:** AL3198HV: 100 assay points AL3198C: 500 assay points AL3198F: 5000 assay points

The number of assay points is based on an assay volume of 50 µL in 96- or 384-well assay plates using components at the recommended concentrations.

Manufacturing date: June 10, 2024 **Document version:** 1**Product Information****Application:** This kit is designed for the quantitative determination of HEK 293 Host Cell Protein (HCP) using a homogeneous no wash AlphaLISA assay.**Kit contents:** The kit contains 5 components: AlphaLISA Acceptor beads coated with anti-HEK 293 HCP polyclonal antibodies (pAb), Streptavidin-coated Donor beads, Biotinylated anti-HEK 293 HCP pAb, Lyophilized HEK 293 HCP and 10X AlphaLISA HiBlock Buffer.**Sensitivity:** Lower Detection Limit (LDL): 0.302 ng/mL
Lower Limit of Quantification (LLOQ): 1.013 ng/mL**Dynamic Range:** 0.302 – 1,000 ng/mL**Storage:** Store kit in the dark at 4 °C. For reconstituted analyte, aliquot and store at -20 °C. Avoid freeze-thaw cycles.**Stability:** This kit is stable for at least 18 months from the date of manufacture when stored in its original packaging and the recommended storage conditions.**Quality Control**

Lot to lot consistency is confirmed in an AlphaLISA assay. Maximum and minimum signals, EC₅₀ and LDL were measured on the EnVision Multilabel Plate Reader with Alpha option. We certify that these results meet our quality release criteria. Maximum counts may vary between bead lots and the instrument used, with no impact on LDL measurement.

EC ₅₀ :	158.200 ng/mL
LDL:	0.349 ng/mL
LLOQ:	1.131 ng/mL
Min counts:	383 counts
Max counts:	30877 counts

The information provided in this document is valid for the specified lot number and date of analysis. This information is for reference purposes only and does not constitute a warranty or guarantee of the product's suitability for any specific use. Revvity, Inc., its subsidiaries, and/or affiliates (collectively, "Revvity") do not assume any liability for any errors or damages arising from the use of this document or the product described herein. REVVITY EXPRESSLY DISCLAIMS ALL WARRANTIES, INCLUDING WARRANTIES OF MERCHANTABILITY OR FITNESS FOR A PARTICULAR PURPOSE, REGARDLESS OF WHETHER ORAL OR WRITTEN, EXPRESS OR IMPLIED, ALLEGEDLY ARISING FROM ANY USAGE OF ANY TRADE OR ANY COURSE OF DEALING, IN CONNECTION WITH THE USE OF INFORMATION CONTAINED HEREIN OR THE PRODUCT ITSELF.

AL3198-R Rev01

The Revvity logo is displayed in a lowercase, sans-serif font.

Revvity, Inc.
940 Winter Street
Waltham, MA 02451 USA

(800) 762-4000 www.revvity.com

For a complete listing of our global offices, visit www.revvity.com
Copyright ©2023, Revvity, Inc. All rights reserved.